

# The Ultimate Challenge.... Reverse Engineering Ourselves Cracking The Code Of The Human Operating System



Over 500 protein kinases and 140 protein phosphatases target more than 500,000 phosphorylation sites in humans to transmit and process the commands from thousands of extracellular mediators to regulate gene expression and other cellular processes. Millions of unique connections between these and other signalling proteins ultimately orchestrate the life and death of cells. Many of these protein interactions have been highly conserved in eukaryotic organisms for over a billion years on this planet. But our knowledge of the composition and architecture of these complex signalling systems is still very rudimentary.

Kinexus Bioinformatics Corporation is pleased to participate in the mapping of signal transduction networks through our unique offering of systems proteomics and bioinformatics services to the academic and industrial biomedical research communities. With the release of this poster, we are proud to announce the launch of PhosphoNET, a free on-line resource from Kinexus at [www.PhosphoNET.ca](http://www.PhosphoNET.ca) with detailed information about tens of thousands of known phosphorylation sites. To familiarize yourself with many known signalling pathways, download our complementary Pathway ScreenSaver from our website at [www.kinexus.ca](http://www.kinexus.ca).

[www.kinexus.ca](http://www.kinexus.ca)

1-866-KINEXUS

[info@kinexus.ca](mailto:info@kinexus.ca)